

# SARTORIUS

## Simplifying Progress

Microsart® ATMP Extraction  
Microsart® ATMP Bacteria & Fungi

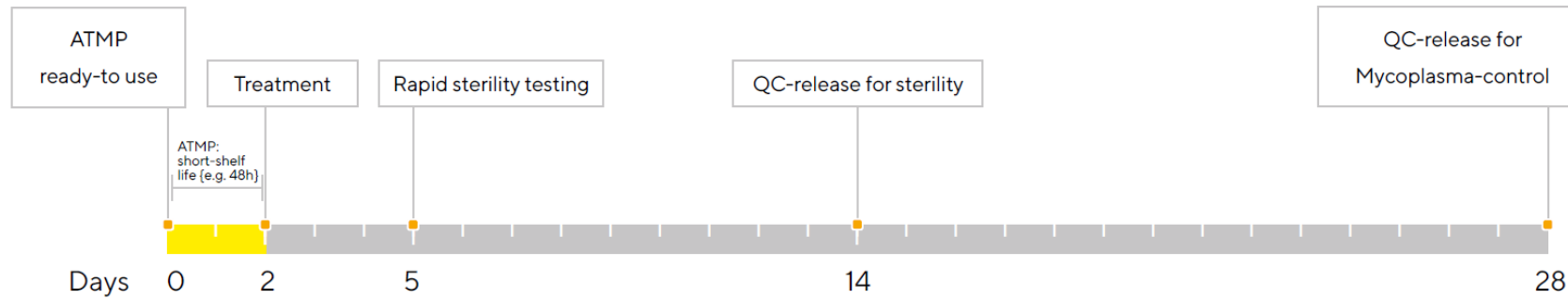
Microsart® ATMP Sterile Release

2022

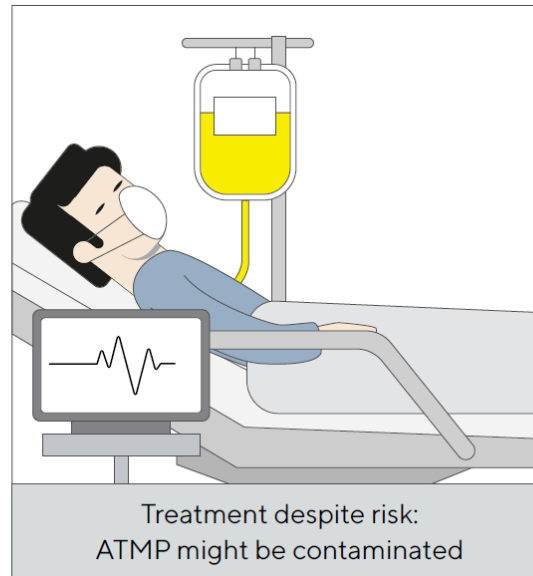
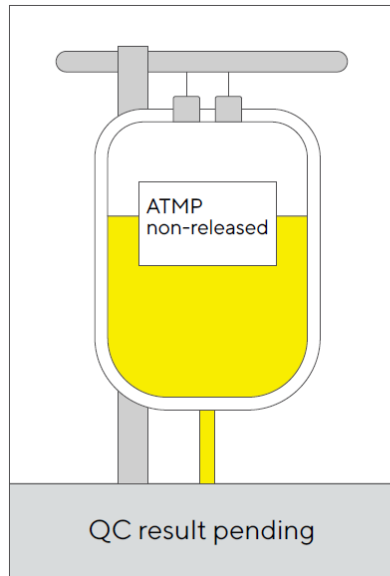


Why new methods?

# ATMPs put microbiological QC to novel challenges

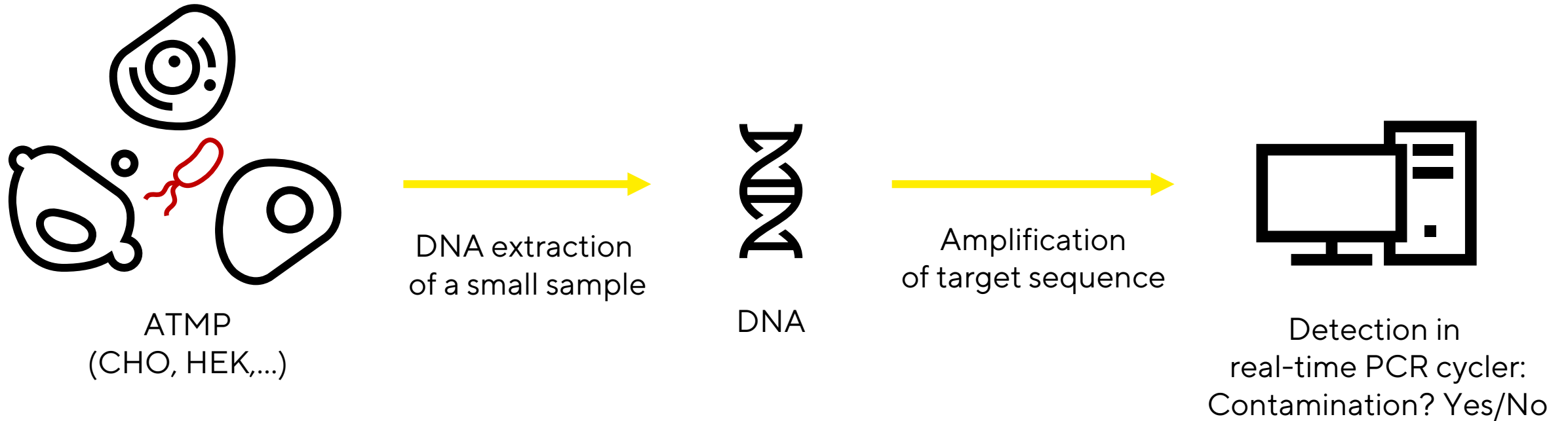


5, 14 or 28 days  
of waiting  
is too long  
for ATMPs!



Why new methods?

## Nucleic acid techniques



**Results within 3 hours!**

# Bacteria & fungi contamination detection

- Real-time PCR allows detection of bacteria and fungi
  - In 3 h
  - Down to 2.5-99 CFU/ml
- Validated combination
  - In accordance with EP 5.1.6, USP 1223, EP 2.6.27, and USP 1071
- Support
  - *Product Validation Report* containing all experimental details
  - *Matrix Validation Proposal* giving an overview of the required set up and materials
  - *Matrix Validation Template* containing detailed information for the customer specific matrix validation
  - **Technical support** during matrix validation process



Microsart® ATMP Extraction



Microsart®  
ATMP Bacteria



Microsart®  
ATMP Fungi



Microsart® ATMP Sterile Release

# Workflow bacteria & fungi contamination detection

- DNA isolation using the Microsart® ATMP Extraction kit
  - Extraction protocol includes centrifugation step to remove free bacterial DNA
  - Harsh extraction allows to isolate Fungi & Bacteria
- 2 real-time PCRs using the Microsart® ATMP Fungi & Microsart® ATMP Bacteria kit
  - Taq-Man® System → reduce false-positive signals
  - Duplex real-time PCR assay → reduce false-negative signals
  - Universal assay for different real-time PCR cycler → FAM™ and ROX™
  - Highly stability & no freezing → Lyophilized reagents



Microsart® ATMP Extraction



Microsart® ATMP Bacteria



Microsart® ATMP Fungi



Microsart® ATMP Sterile Release



Simplifying Progress

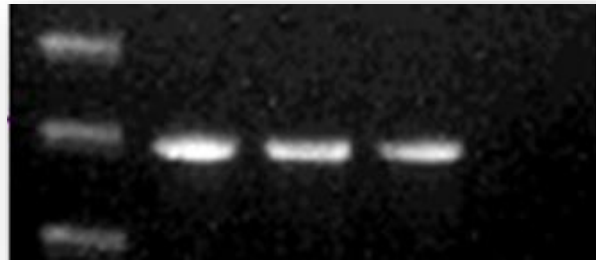
Technical background  
DNA-based detection methods

**SARTORIUS**

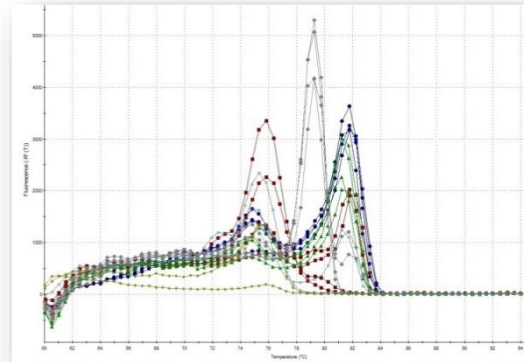
# DNA-based detection methods

real-time PCR

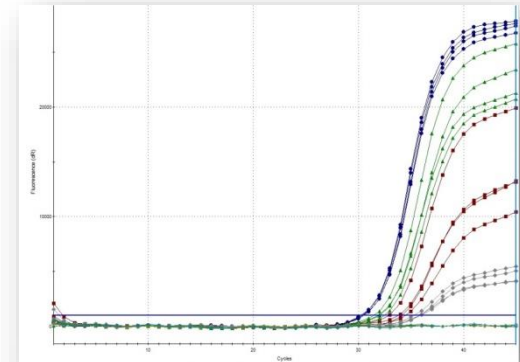
Conventional PCR



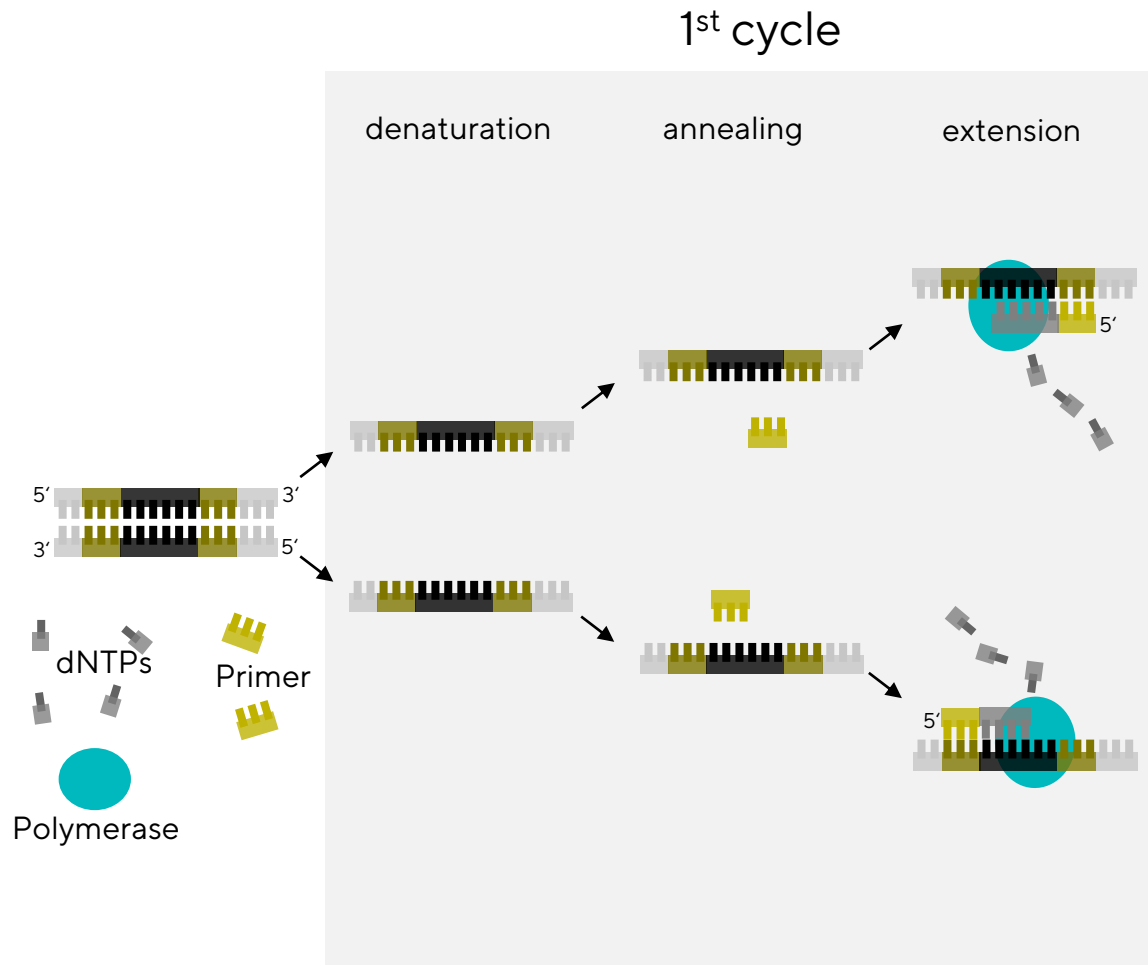
SYBR Green I



TaqMan® Probe

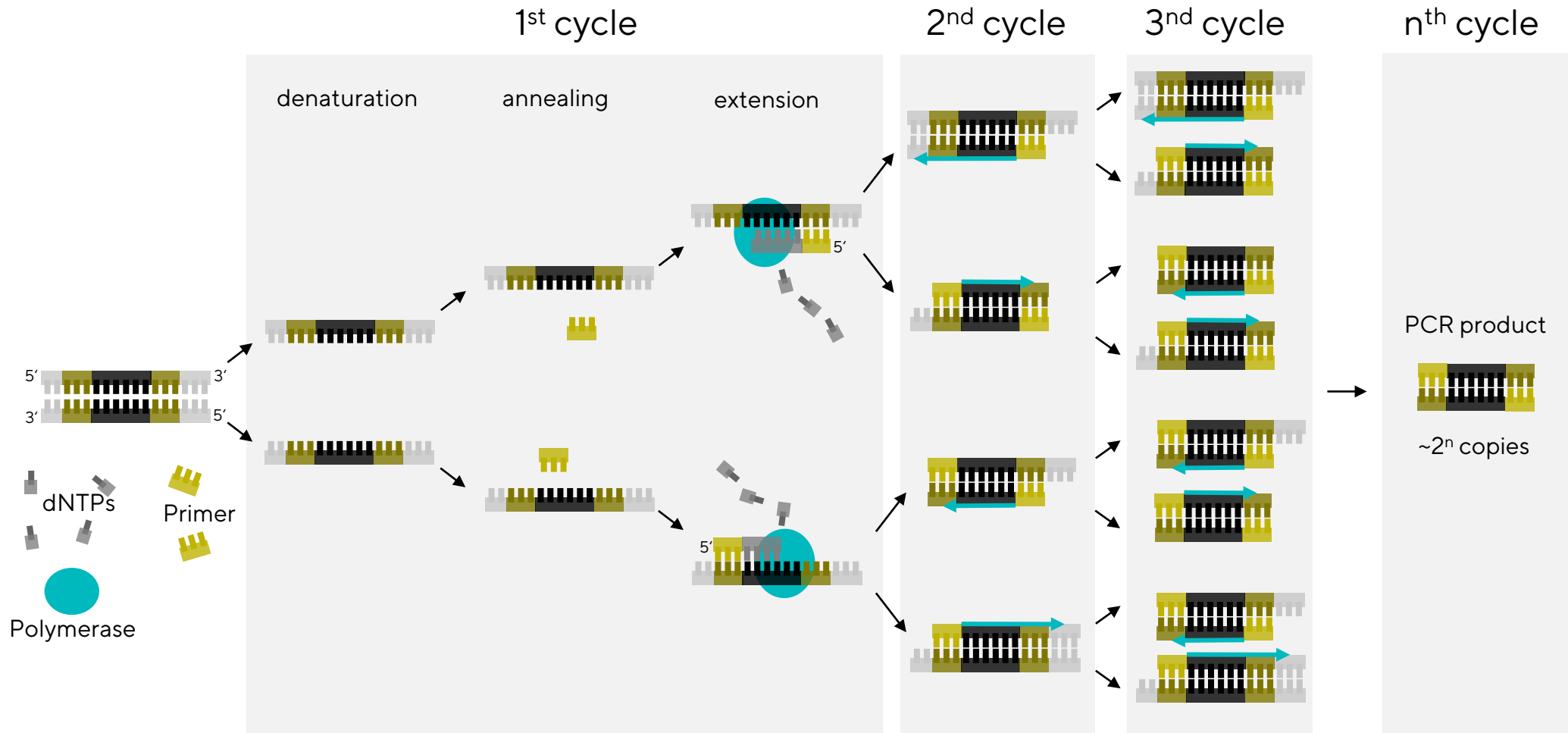


# What is a conventional PCR?

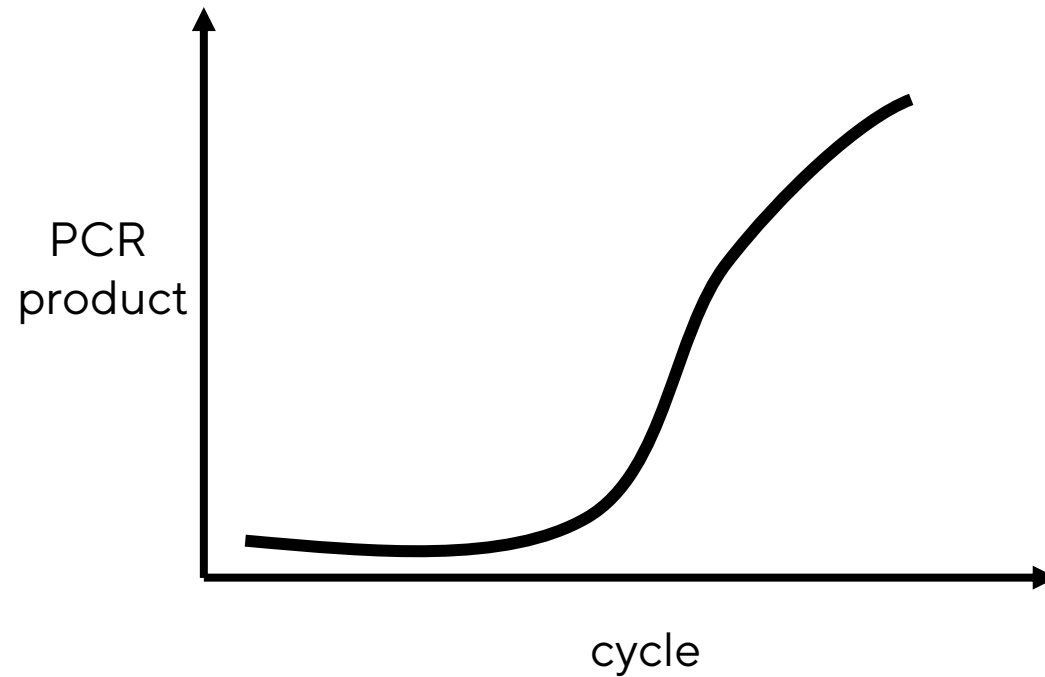




# What is a conventional PCR?

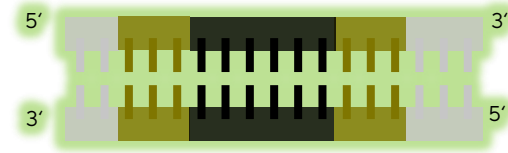
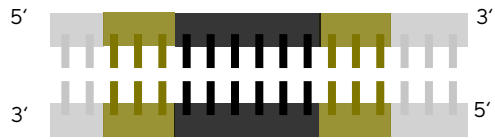
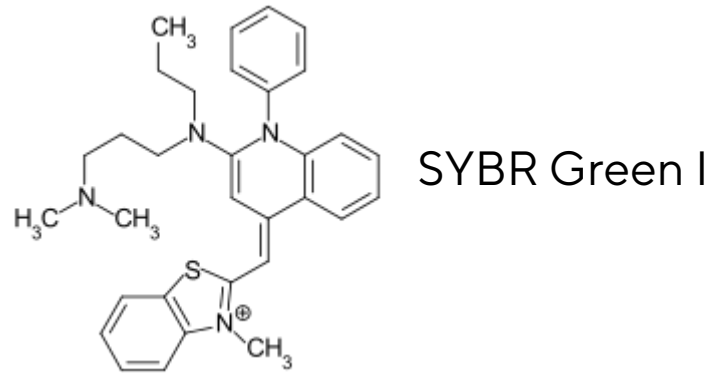


# A real-time PCR visualizes the reaction

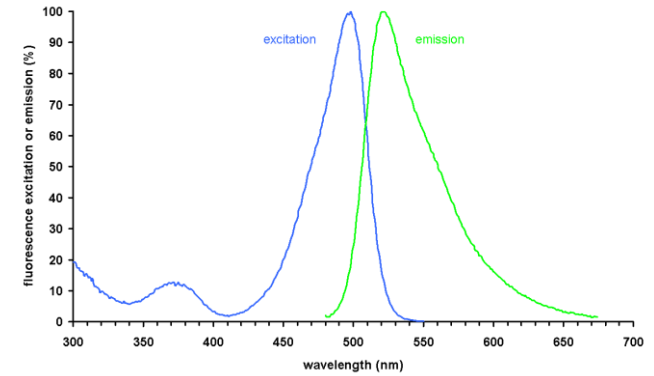


How does that work?

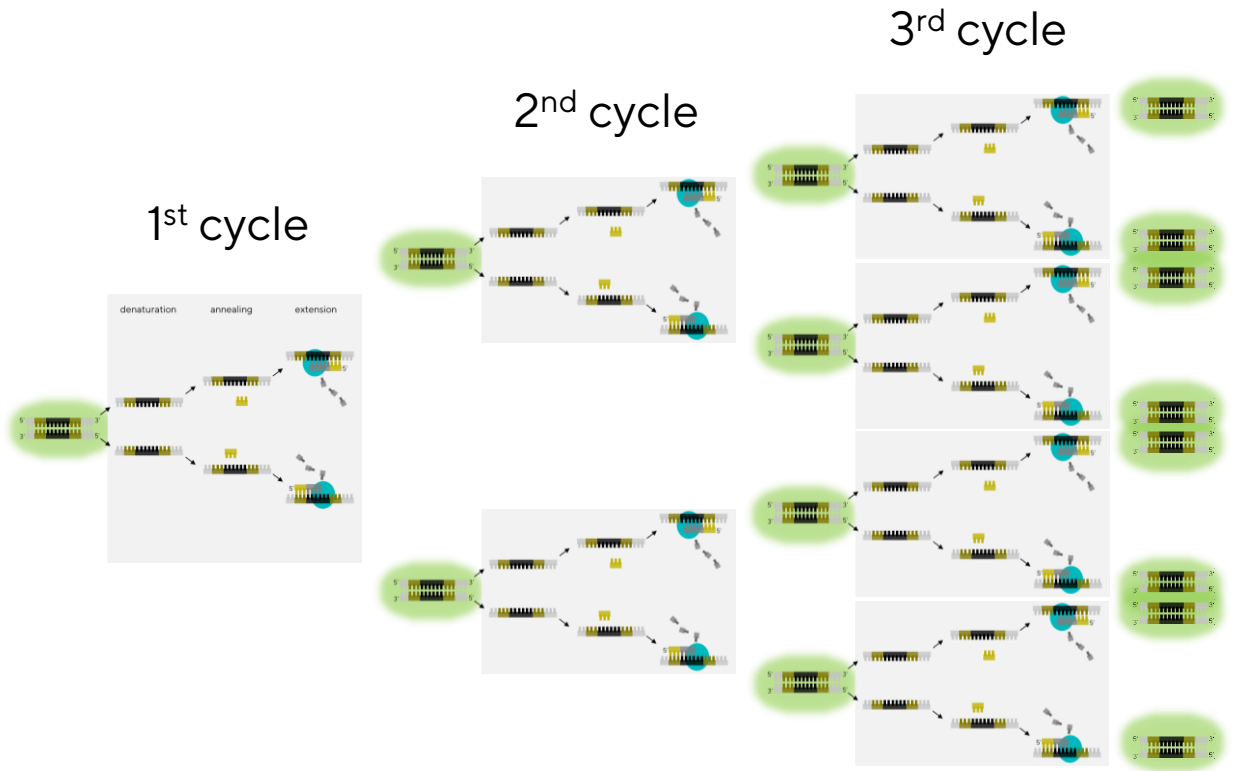
# Real-time PCR



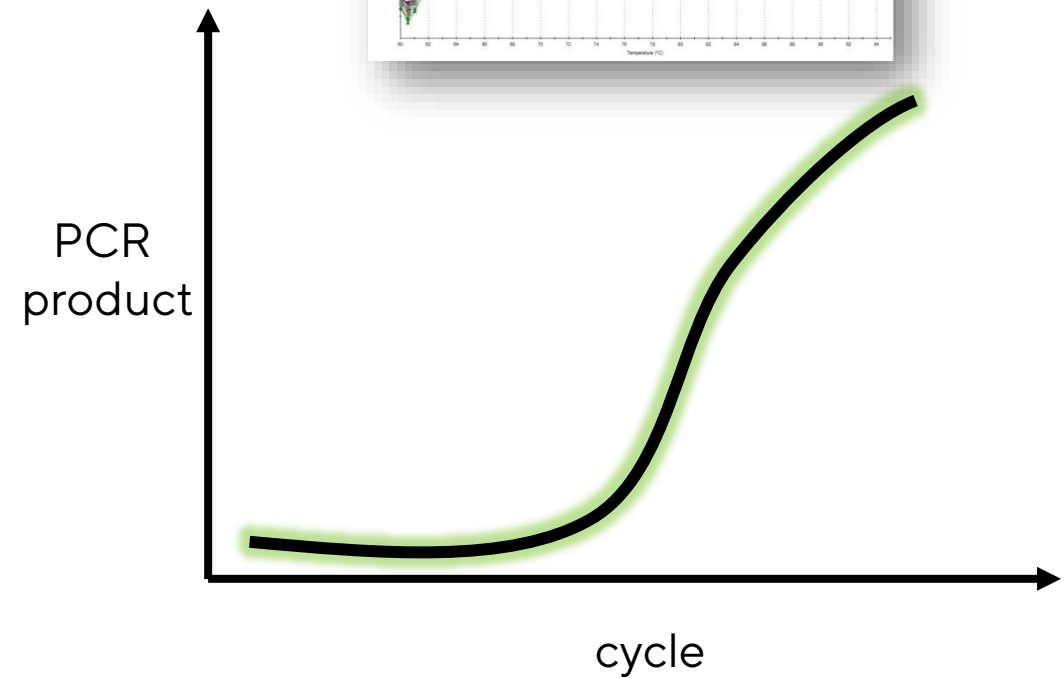
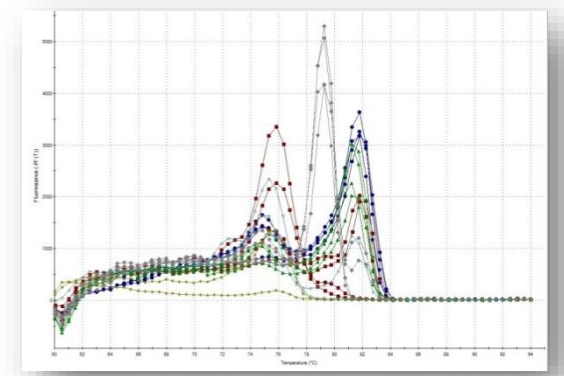
The dye SYBR Green I binds to double stranded DNA!



# Real-time PCR

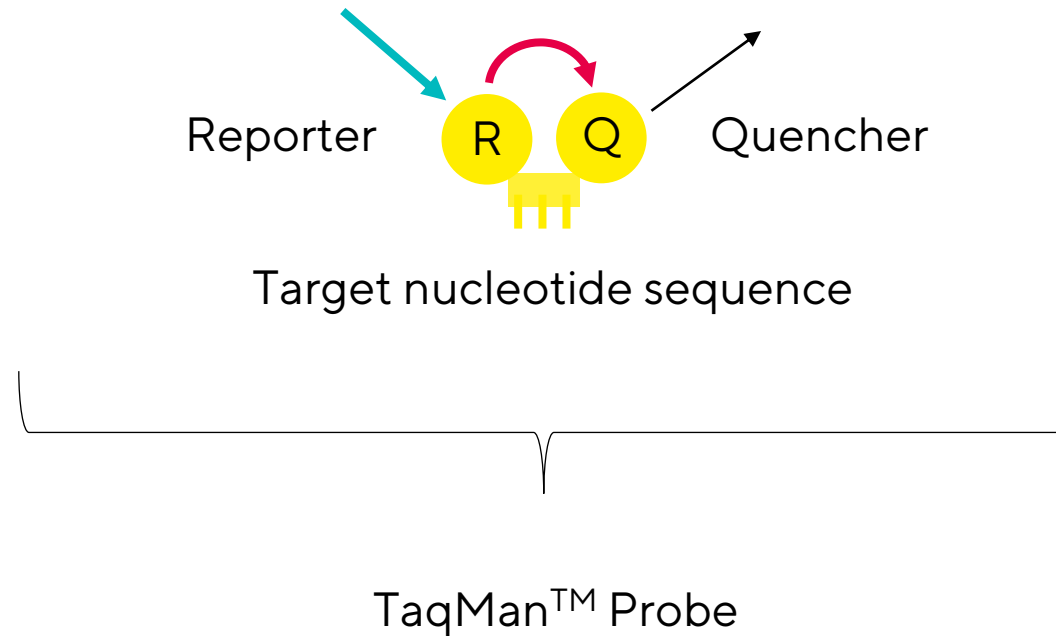


SYBR Green I



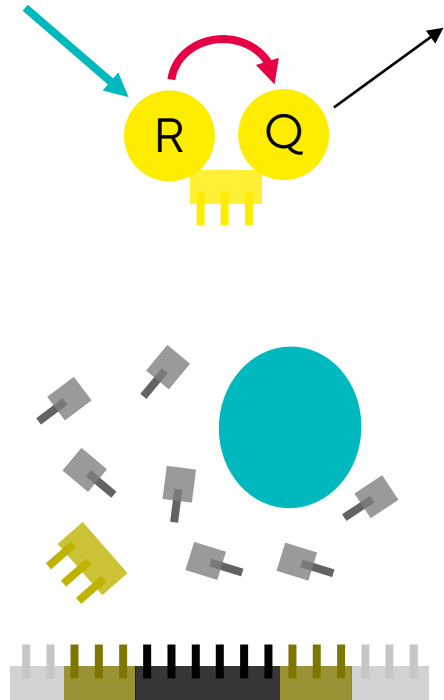
Unspecific binding of SYBR Green I can result in false-positive signals!

# A TaqMan™ probe is more specific compared to SYBR Green I

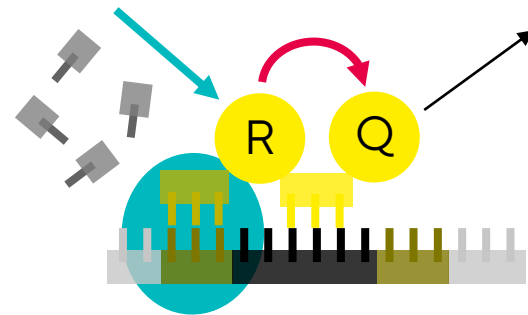


# TaqMan™ real-time PCR

TaqMan™ probe is degraded during real-time PCR

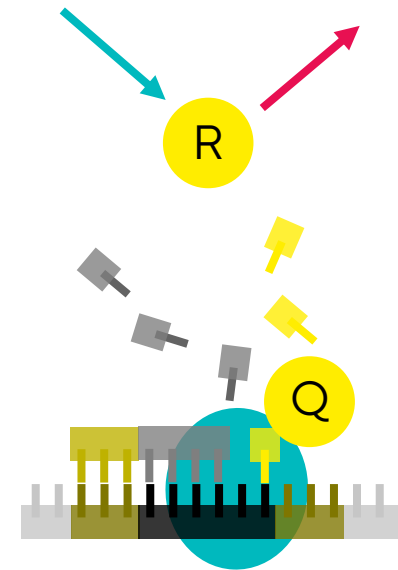


As long as the probe is complete no light signal can be detected



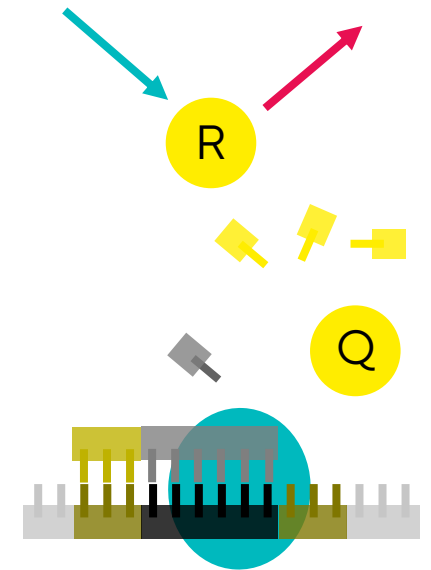
Taq Polymerase functions:

- DNA amplification
- 5'-3' exonuclease activity

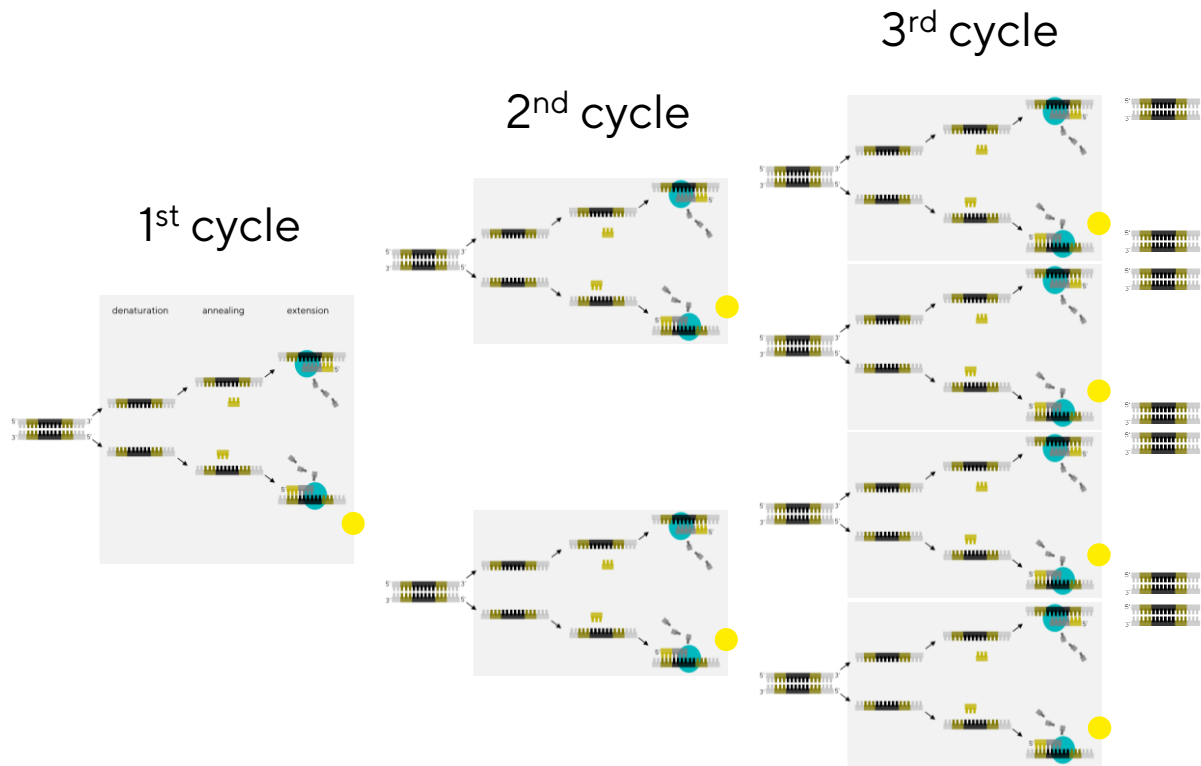


During elongation:

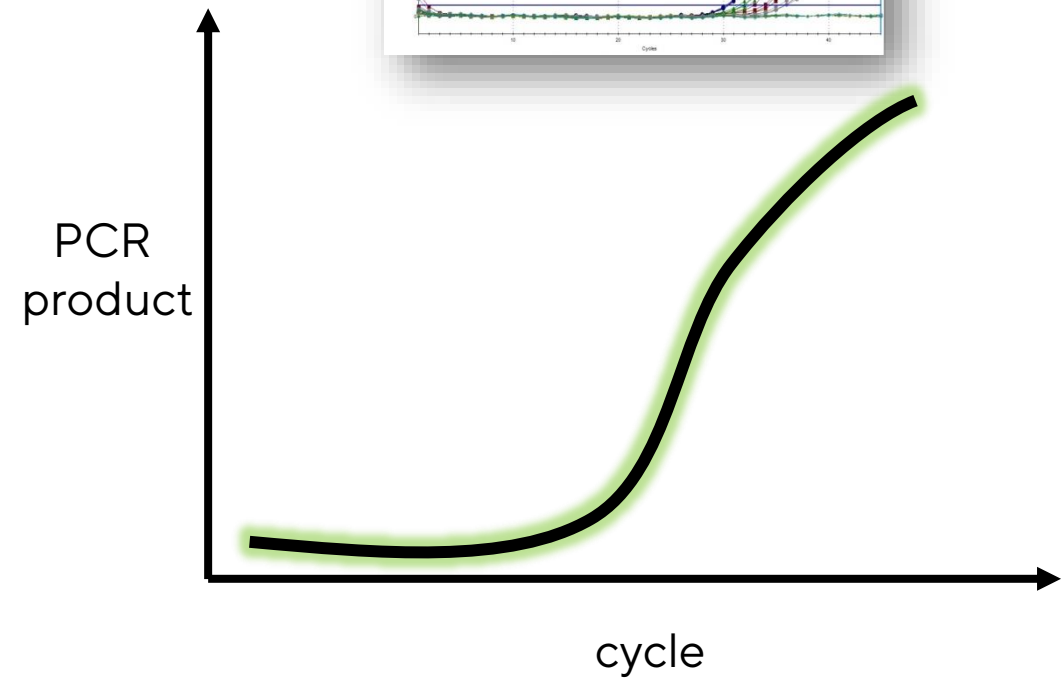
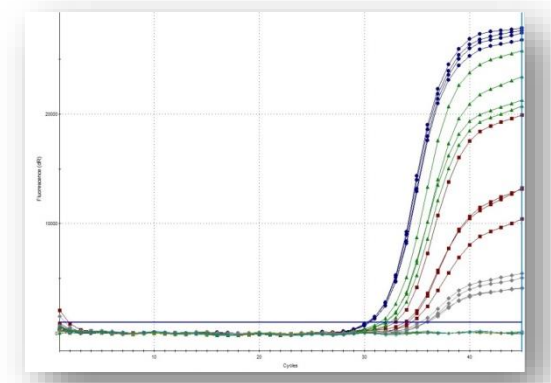
- Polymerase hydrolyses probe
- Dye and quencher are separated
- Reporter dye emits light signal



# TaqMan™ real-time PCR



TaqMan® Probe



The specificity of TaqMan™ system **reduces false-positive results!**

# A duplex real-time PCR assay monitors PCR functionality

Problem:

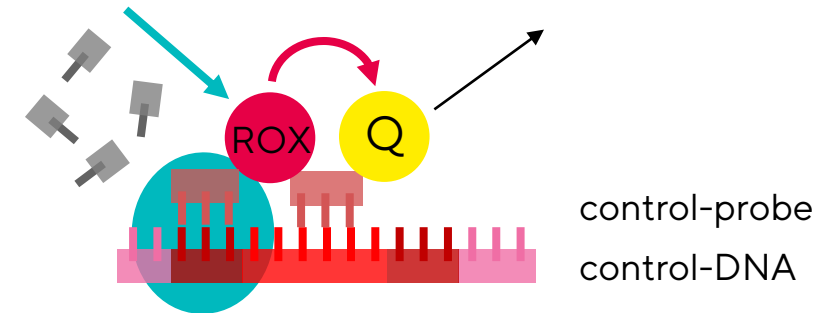
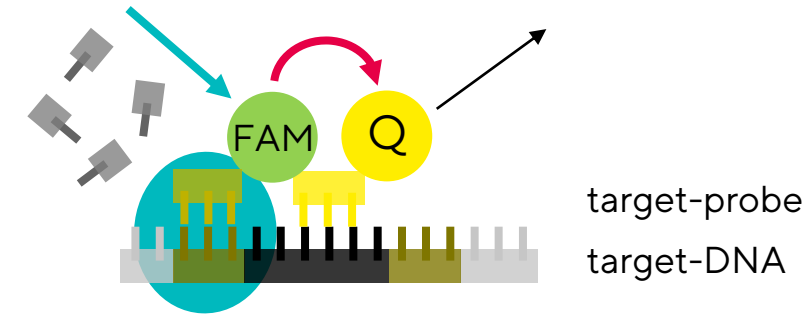
What does **no signal** mean?  
No microbial contamination?  
PCR inhibition?

Solution:

Include a second real-time PCR and  
a control DNA that must lead to a  
signal!

→ If this internal control reaction  
does not lead to a signal,  
the PCR is inhibited.

**Duplex assay**  
= two independent real-time PCRs  
in one run using different  
fluorophores



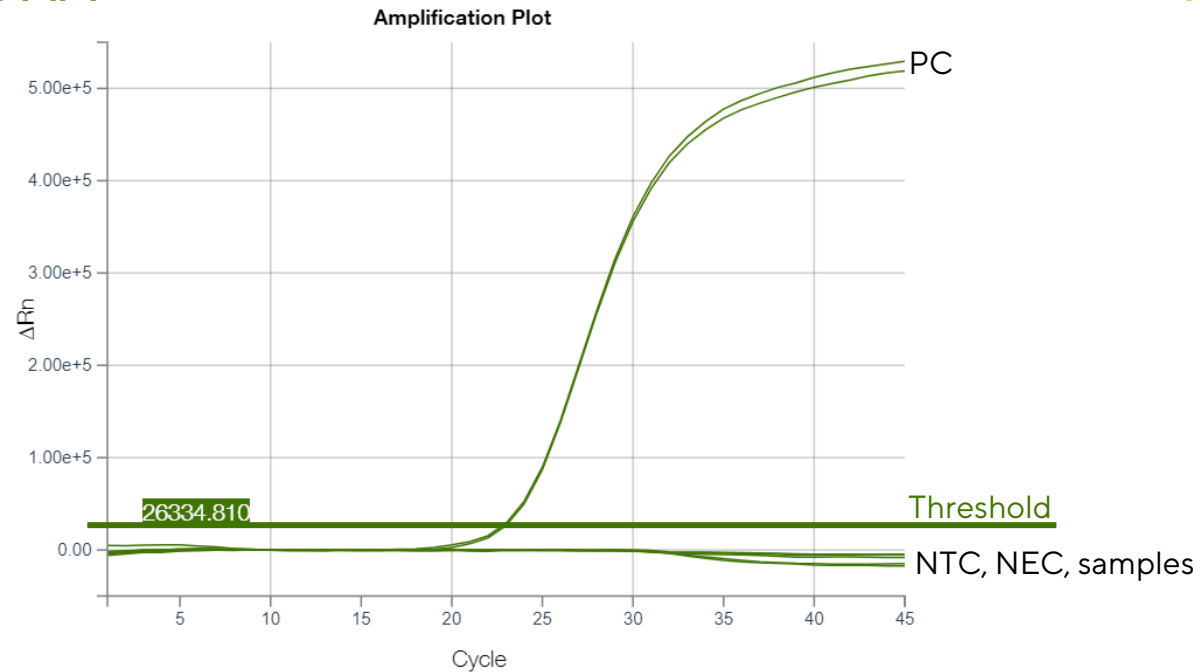
The internal control reaction **reduces false-negative results!**



# A duplex real-time PCR Analysis

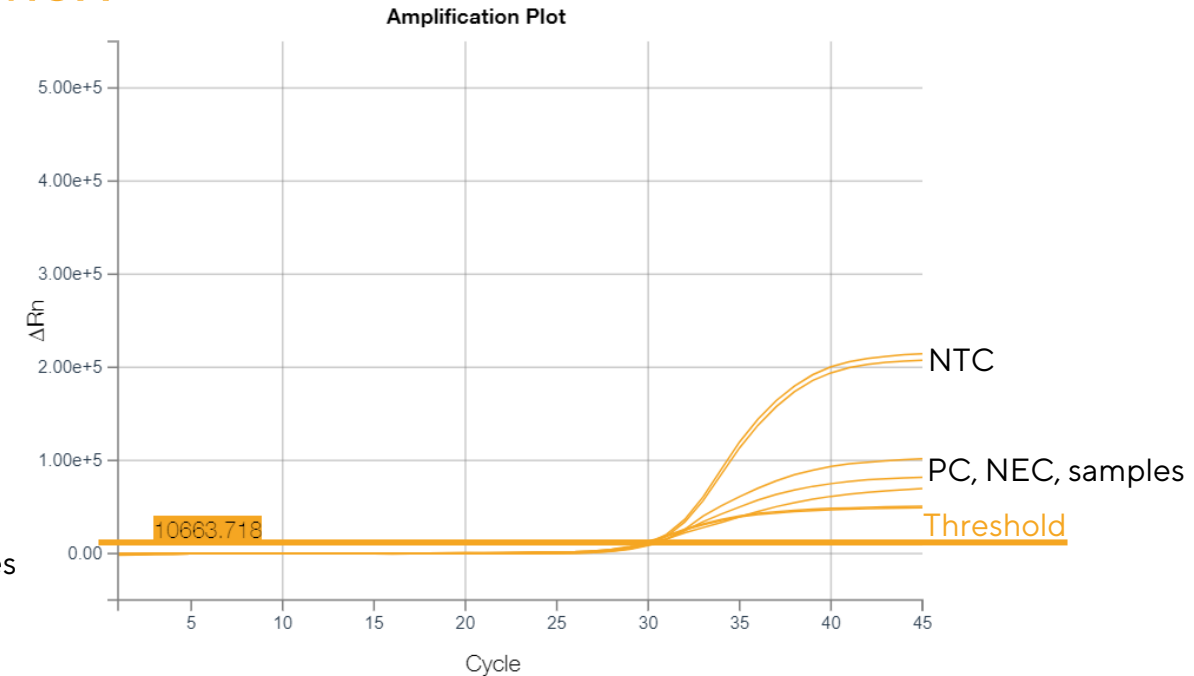
This analysis is done for both, the bacteria real-time PCR and the fungi real-time PCR

FAM



There is no contamination in the samples, because only the positive control is detectable in the FAM channel

ROX



There is no PCR inhibition, because the internal control DNA was detected in all reactions.

## What is in the kits?

### real-time PCR master mix Bacteria

- Primer for bacterial DNA
- Primer for control DNA
- FAM probe for target DNA
- ROX probe for control DNA
- Taq polymerase
- Buffer

### real-time PCR master mix Fungi

- Primer for fungal DNA
- Primer for control DNA
- FAM probe for target DNA
- ROX probe for control DNA
- Taq polymerase
- Buffer



Rehydration buffer

Internal Control DNA

Positive Control DNA

Ultrapure Water

Elution Buffer\*

Lysis Buffer\*

\*Only in Microsart® ATMP Sterile Release



Simplifying Progress




Hands on fungi & bacteria detection


**SARTORIUS**


# Microsart® ATMP Extraction



## 1. Sample Collection



 + 1 ml  
sample material  
Processing tube




 15 min  
≥ 16,200 × g

  
Discard  
supernatant



Store at ≤ -18 °C  
  
or proceed to  
DNA extraction

## 2. DNA Extraction

 + 500 µl  
Lysis Buffer  
(transparent cap)  


 ≥ 30 sec vigorously  
 80 °C, 10 min  
 ≥ 16,200 × g, 10 min

  
Remove supernatant  
carefully

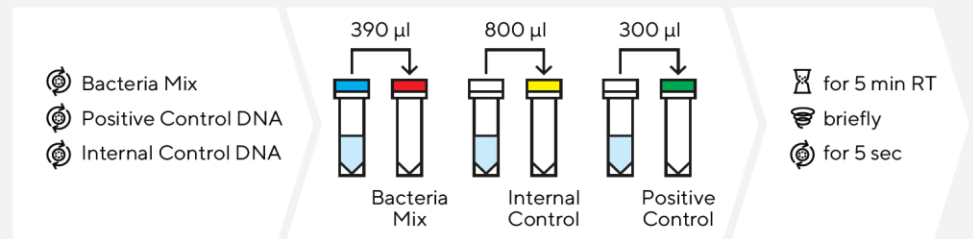
 + 100 µl Suspension  
Buffer (violet cap)  
 ≥ 30 sec vigorously

  
DNA ready for PCR

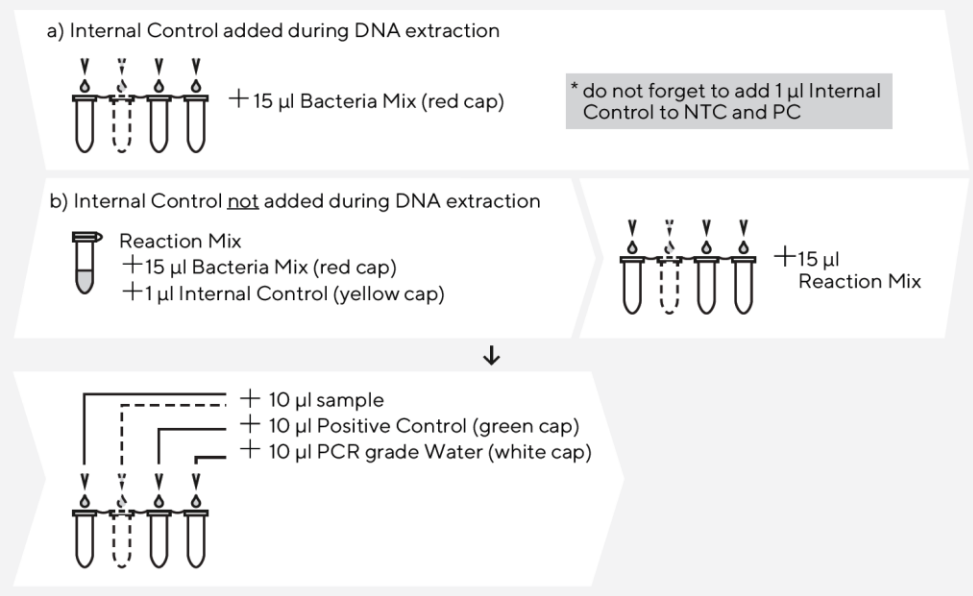
# Microsart® ATMP Bacteria



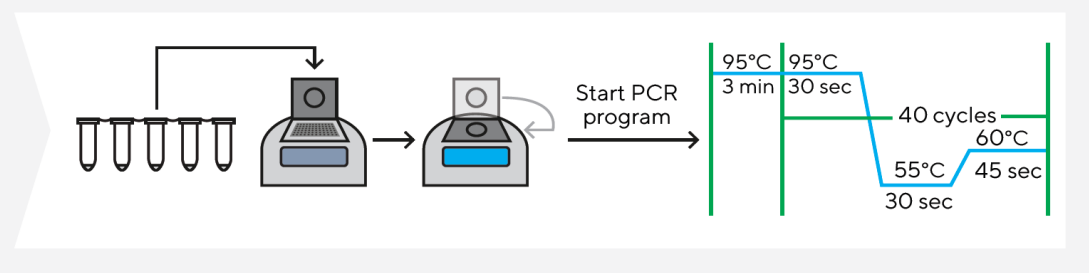
## 1. Rehydration of Reagents



## 2. Preparation of PCR Reaction



## 3. Starting PCR Reaction



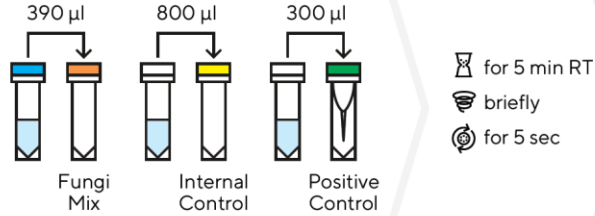
- Rehydration Buffer  
■ Bacteria Mix  
 PCR grade Water  
■ Positive Control  
■ Internal Control
- ⌚ incubate  
🌀 vortex  
🌀 centrifuge  
+ add
- storage +2 - +8 °C  
 after rehydration ≤ -18 °C

This procedure overview is not a substitute for the detailed manual. ST\_SI\_Microsart®-ATMP-Bacteria\_04\_EN

# Microsart® ATMP Fungi

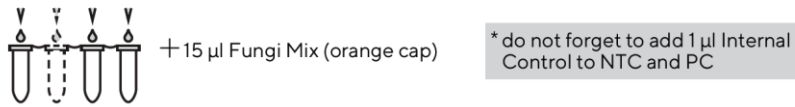
## 1. Rehydration of Reagents

- 🌀 Fungi Mix
- 🌀 Positive Control DNA
- 🌀 Internal Control DNA

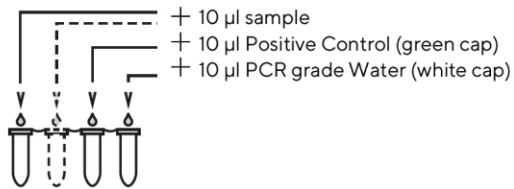


## 2. Preparation of PCR Reaction

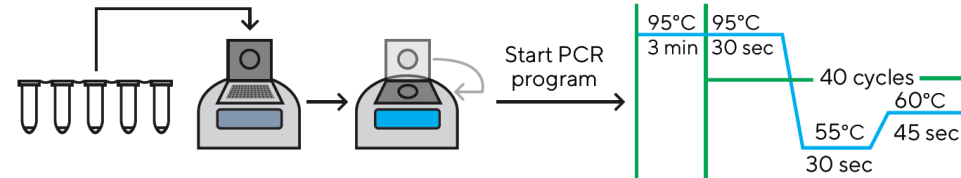
### a) Internal Control added during DNA extraction



### b) Internal Control not added during DNA extraction



## 3. Starting PCR Reaction



- Rehydration Buffer
- Fungi Mix
- PCR grade Water
- Positive Control
- Internal Control

- ⌚ incubate
- 🌀 vortex
- 🌀 centrifuge
- ⊕ add

storage +2 - +8 °C  
after rehydration ≤ -18 °C

This procedure overview is not a substitute for the detailed manual.

ST\_SI\_Microsart®-ATMP-Fungi\_O2\_EN

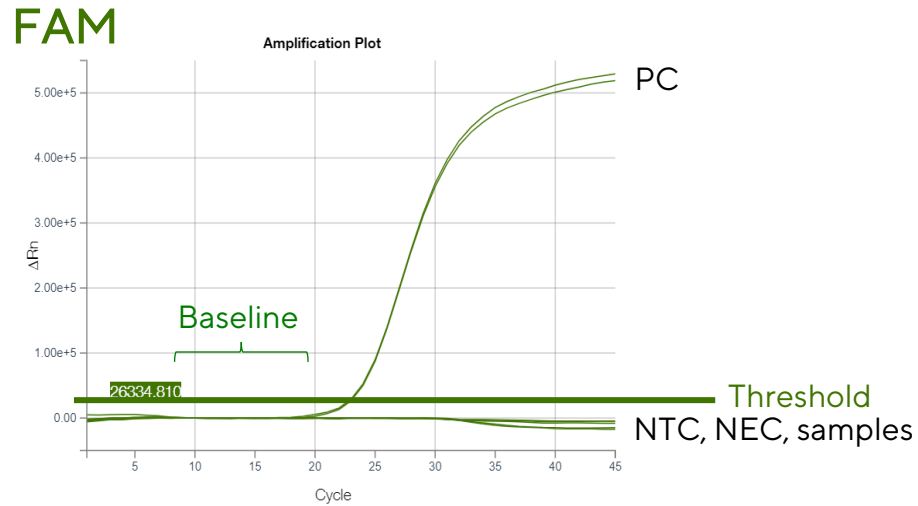


Fungi Mix

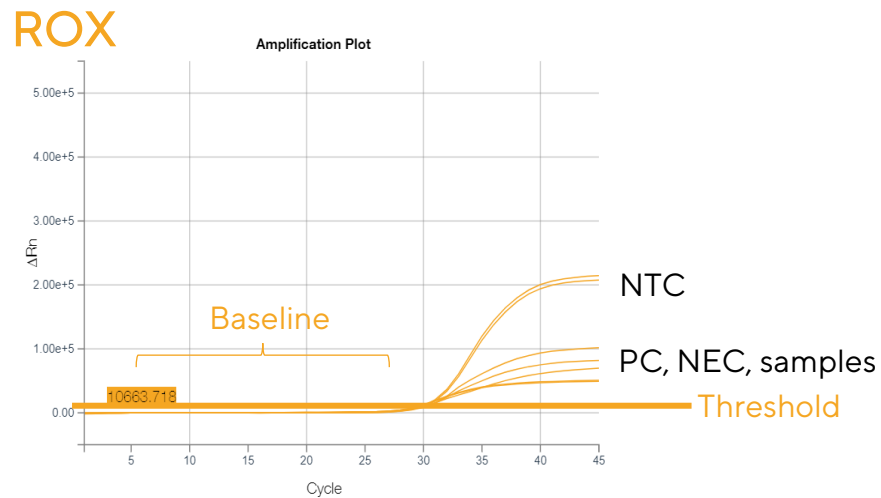
All identical, but the fungi master mix has an orange cap!



# Microsart® ATMP Sterile Release - Analysis



1. Set the baseline to level the curves
2. Set the threshold
3. Check if all controls are as expected  
→ see next slide
4. Analyze your samples











Identical analysis workflow for both, the fungi and the bacteria real-time PCR!





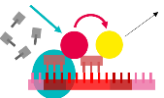

# Result interpretation




The result interpretation is identical for both, the bacteria real-time PCR and the fungi real-time PCR

Templates added to the real-time PCR

	PC Positive Control	NTC No Template Control	Sample	NEC Negative Extraction Control
FAM Template				
ROX Template				

Read-outs

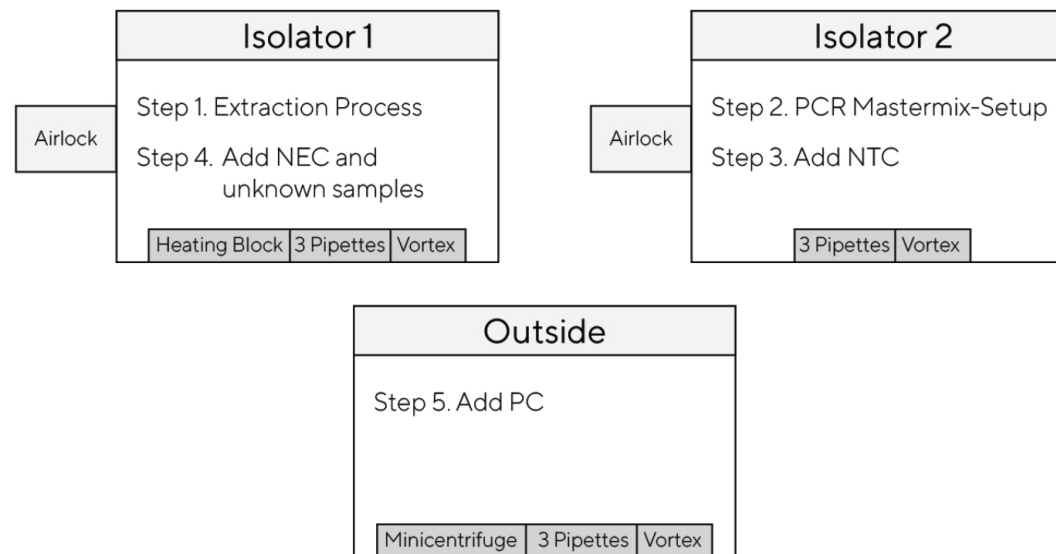
Target	 POSITIVE	 NEGATIVE	To be tested!	 NEGATIVE
Internal Control	 Does not matter	 POSITIVE		 POSITIVE

 PCR grade Water  
 Positive Control  
 Internal Control

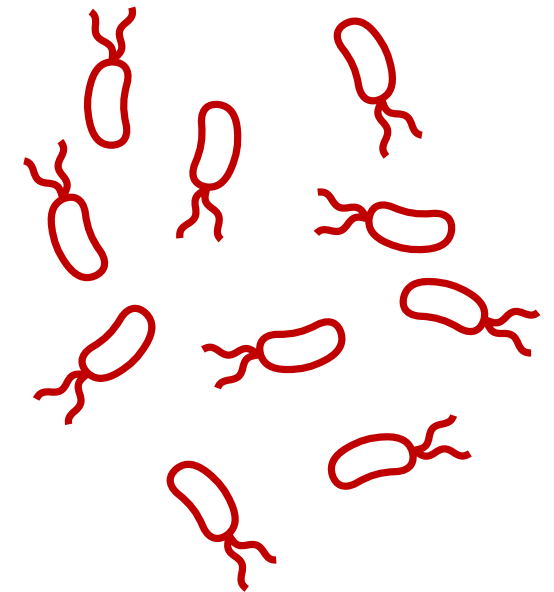


# Tips and tricks

- Extensive cleaning with chlorine-based agents
- Avoid cleaning with ethanol → Ethanol precipitates DNA
- Spatial separation of DNA extraction process, master mix setup and positive control
  - E.g. isolator/glovebox or laminar flow
- Work carefully e.g. do not touch the lids of open tubes



**Bacterial DNA  
is everywhere!**



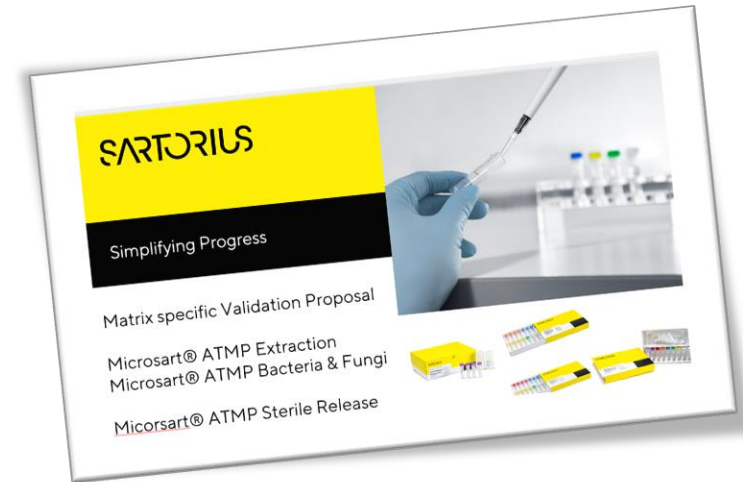
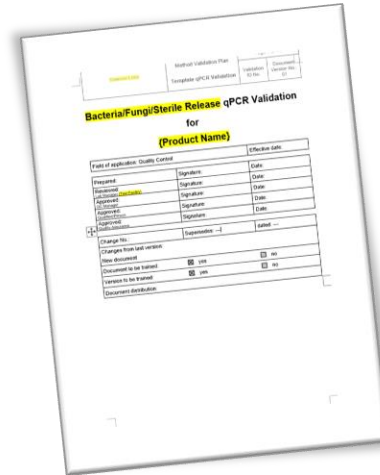
Simplifying Progress



Sartorius validation support

**SARTORIUS**

# Validation reports, templates & testing proposals



- Product Validation Reports
  - Microsart® ATMP Bacteria + Microsart® ATMP Extraction
  - Microsart® ATMP Fungi + Microsart® ATMP Extraction

- Validation Template
  - Combined validation Template for Microsart® ATMP Bacteria + Microsart® ATMP Fungi + Microsart® ATMP Extraction

- Validation proposal
  - Standard matrix specific validation
  - Individual support

## Further support for your validation

- Microsart® Validation Standard (99 CFU/Vial) & Microsart® Calibration Reagents (10<sup>8</sup> GC/Vial for bacteria, 10<sup>6</sup> GC/Vial for fungi)

- *Bacillus subtilis*
- *Pseudomonas aeruginosa*
- *Kocuria rhizophila* | *Micrococcus luteus*
- *Clostridium sporogenes*
- *Bacteroides vulgatus*
- *Staphylococcus aureus*
- *Candida albicans*
- *Aspergillus brasiliensis*
- *Aspergillus fumigatus*
- *Penecillium chrysogenum*
- *Candida glabrata*
- *Candida krusei*
- *Candida tropicalis*

Non-viable CFU Standards!

DNA Standards!

Do you miss a species? Let us know!

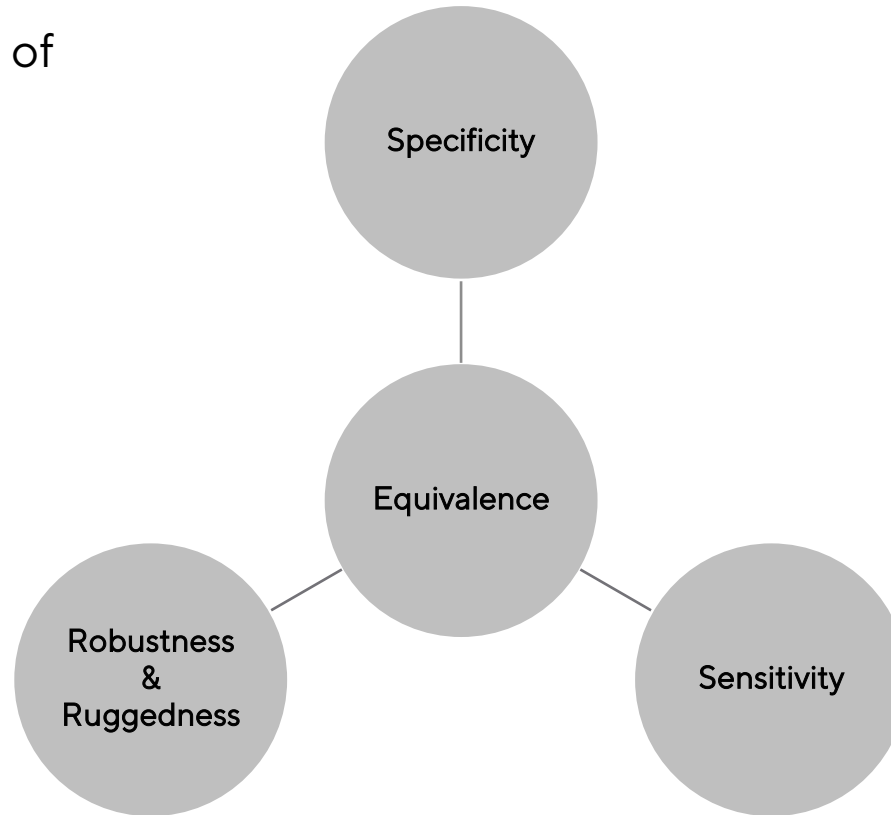
# Status quo regulatory landscape

Microbiological QC-Release testing			
Method	Mycoplasma	Bacteria	Fungi
Classical testing	USP<63>   EP 2.6.7 28 days	USP<71>   EP 2.6.1 Sterility testing 14 days	
real-time PCR-based	EP 2.6.7 (USP<1223>/EP 5.1.6)	USP<1071> EP 2.6.27 (USP<1223>/EP 5.1.6)	

# Validation overview

Regulatory guidance for validation of rapid / alternativ methods:

- PDA, TR 33
- USP <1223>
- USP <1071>
- EP 5.1.6 part 4-1-1 primary validation by supplier
- EP 2.6.27
- (USP<71>)
- (EP2.6.1)



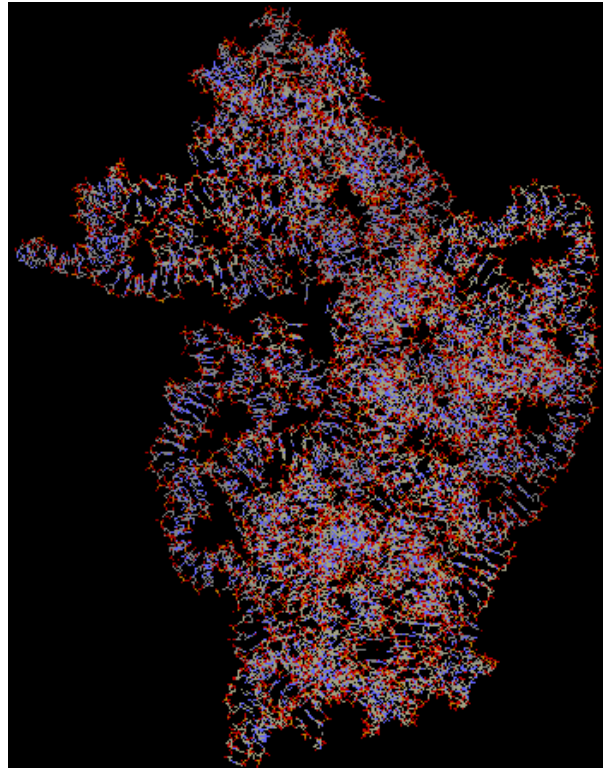
NOT a replacement for the classical sterility test,  
but valuable addition for patient safety

Including Guidance of the  
German Governmental  
Regulatory Agency (part  
of EMEA)

Paul-Ehrlich-Institut 

# *In silico* prediction by sequence alignment and blast

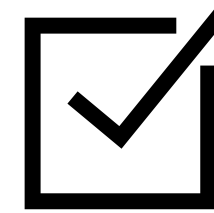
16S / 18S rDNA



<http://www.biochem.umd.edu/biochem/kahn/bchm465-01/ribosome/16SrRNA.html>



<https://www.ncbi.nlm.nih.gov/>



Specificity

## Detection range – Microsart® ATMP Bacteria

Species	Strain No	Primer Mismatches		
		Forward primer	Probe	Reverse primer
<i>Bacillus subtilis</i>	ATCC 6633	0	0	0
<i>Clostridium sporogenes</i>	ATCC 3584	0	0	0
<i>Pseudomonas aeruginosa</i>	ATCC 9027	0	0	0
<i>Staphylococcus aureus</i>	ATCC 6538	0	0	0

	Primer Mismatches			
	0	1	2	3
Bacteria	48.8 %	69.4 %	85.7 %	94.7 %
Archaea	n.a	n.a	0.1%	40.4 %
Eukaryotes	0%	0%	0.1%	0.3 %

Accepting 3 primer mismatches, **94.7 % of the bacteria** are detected





# Detection range – Microsart® ATMP Fungi

Accepting 2 primer mismatches, already **37 % of the fungi** are detected, including all species of clinical and bioprocess relevance.



Genus	Coverage
<i>Alternaria</i>	97.7 %
<i>Aspergillus</i>	95.3 %
<i>Aureobasidium</i>	93.5 %
<i>Bipolaris</i>	98 %
<i>Candida</i>	86.3 %
<i>Chaetomium</i>	3.6%
<i>Cladosporium</i>	95.5 %
<i>Curvularia</i>	100 %
<i>Epidermophyton</i>	100 %
<i>Exserohilum</i>	97.4 %
<i>Fusarium</i>	95.9 %
<i>Memnoniella (Stachybotrys)</i>	86,7 %
<i>Microsporum</i>	100 %
<i>Myrothecium</i>	100 %
<i>Paecilomyces</i>	100%
<i>Penicillium</i>	98.2 %
<i>Malassezia*</i>	0.1 %
<i>Rhizopus</i>	4 %
<i>Scopulariopsis</i>	0 %
<i>Trichoderma</i>	98 %
<i>Trichophyton</i>	100 %

← soil, air, plant debris

← skin microbiome  
 ← organic substances  
 ← soil, decaying wood

Species	Strain No	Primer Mismatches		
		Forward primer	Probe	Reverse primer
<i>Aspergillus brasiliensis</i>	ATCC 6275	0	0	0
<i>Candida albicans</i>	ATCC 18804	0	0	0



# Matrix Effects

## Bacteria

		Results
Hela	No C <sub>q</sub>	0/2
	No C <sub>q</sub>	
Vero	No C <sub>q</sub>	0/2
	No C <sub>q</sub>	
CHO-K1	No C <sub>q</sub>	0/2
	No C <sub>q</sub>	
RK13	No C <sub>q</sub>	0/2
	No C <sub>q</sub>	
CHO-DG44	No C <sub>q</sub>	0/2
	No C <sub>q</sub>	
CHO XM111-10	No C <sub>q</sub>	0/2
	No C <sub>q</sub>	
L-9296 (NCTC)	No C <sub>q</sub>	0/2
	No C <sub>q</sub>	

## Fungi

			Results
Hela	No C <sub>q</sub>	No C <sub>q</sub>	0/8
	No C <sub>q</sub>	No C <sub>q</sub>	
	No C <sub>q</sub>	No C <sub>q</sub>	
	No C <sub>q</sub>	No C <sub>q</sub>	
Vero	No C <sub>q</sub>	No C <sub>q</sub>	0/8
	No C <sub>q</sub>	No C <sub>q</sub>	
	No C <sub>q</sub>	No C <sub>q</sub>	
	No C <sub>q</sub>	No C <sub>q</sub>	
CHO-K1	No C <sub>q</sub>	No C <sub>q</sub>	0/8
	No C <sub>q</sub>	No C <sub>q</sub>	
	No C <sub>q</sub>	No C <sub>q</sub>	
	No C <sub>q</sub>	No C <sub>q</sub>	
HPBMC	No C <sub>q</sub>	No C <sub>q</sub>	0/8
	No C <sub>q</sub>	No C <sub>q</sub>	
	No C <sub>q</sub>	No C <sub>q</sub>	
	No C <sub>q</sub>	No C <sub>q</sub>	
Jurkat	No C <sub>q</sub>	No C <sub>q</sub>	0/8
	No C <sub>q</sub>	No C <sub>q</sub>	
	No C <sub>q</sub>	No C <sub>q</sub>	
	No C <sub>q</sub>	No C <sub>q</sub>	

For the tested matrices, **no matrix effects** were detected.

# Limit of detection – Microsart® ATMP Bacteria/Fungi

EP 2.6.1  
USP<71>

regulatory advice;  
EP 2.6.27; user feedback

Currently tested = 6



*Bacillus subtilis*  
*Clostridium sporogenes*  
*Pseudomonas aeruginosa*  
*Staphylococcus aureus*

*Candida albicans*  
*Aspergillus brasiliensis*

23/24 positive

Recommended Extension = 20

+ Colony Forming Units (CFU)

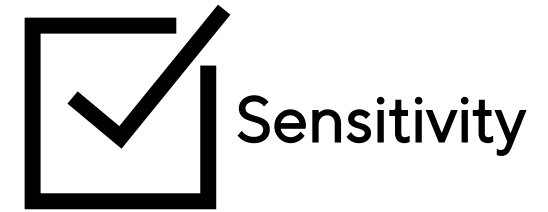
- Streptococcus pyogenes*
- Bacteroides vulgatus*
- Escherichia coli*
- Pseudomonas protegens*
- Bacillus cereus*
- Enterococcus faecalis*
- Kocuria rhizophila*
- Staphylococcus epidermidis*
- Serratia marcescens*
- Cutibacterium acnes*

Genome Copies (GC)

- Candida tropicalis*
- Candida glabrata*
- Candida krusei*
- Aspergillus fumigatus*
- Penicillium chrysogenum*
- Bacteroides fragilis*
- Enterobacter cloacae*
- Klebsiella pneumoniae*
- Clostridium perfringens*
- Yersinia enterocolitica*

8/8 positive

Limit of Detection (LOD<sub>95</sub>): 99, 50, 25, 10, 5, 2.5, 1.25 CFU/ml



# Limit of detection

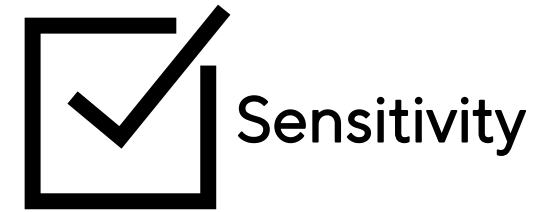
Example: *Candida albicans*

CFU/ml						Mean	
99	Run 1	31.32	31.90	32.78	32.80	32.20	24/24
	Run 2	32.96	32.88	33.07	32.25	32.79	
	Run 3	32.29	32.81	36.03	32.09	33.30	
	Run 4	32.46	32.14	32.29	32.33	32.30	
	Run 5	32.32	32.55	32.71	32.36	34.48	
	Run 6	32.58	31.48	31.89	32.43	32.09	
50	Run 1	33.15	33.79	34.11	34.44	33.87	23/24
	Run 2	34.25	33.86	33.04	35.03	34.04	
	Run 3	33.59	33.90	32.63	33.89	33.50	
	Run 4	34.04	32.94	34.27	33.81	33.76	
	Run 5	No Cq	34.77	32.82	34.48	34.02	
	Run 6	32.58	32.57	32.82	34.81	31.19	
25	Run 1	33.42	34.72	34.88	35.32	34.58	21/24
	Run 2	34.67	37.19	33.28	36.08	35.30	
	Run 3	34.96	37.20	38.18	No Cq	36.78	
	Run 4	37.96	34.90	35.11	34.59	35.64	
	Run 5	33.55	No Cq	35.77	No Cq	34.66	
	Run 6	34.51	35.33	36.17	34.90	35.23	
10	Run 1	35.85	36.48	37.27	37.38	36.74	14/24
	Run 2	No Cq	No Cq	37.00	34.65	35.82	
	Run 3	No Cq	No Cq	No Cq	No Cq	No Cq	
	Run 4	No Cq	35.93	No Cq	37.11	36.52	
	Run 5	35.04	No Cq	No Cq	38.91	36.97	
	Run 6	35.54	36.26	35.84	37.52	36.29	

currently tested

recommended extension

Species	Strain No	Acceptance criterion	LOD <sub>95</sub> (CFU/mL)
<i>Candida albicans</i>	ATCC 10231	23/24	50
<i>Aspergillus brasiliensis</i>	ATCC 16404	23/24	50
<i>Candida tropicalis</i>	ATCC 750	8/8	10
<i>Candida glabrata</i>	ATCC 90030	8/8	25
<i>Candida krusei</i>	ATCC 6258	8/8	50
<i>Aspergillus fumigatus</i>	ATCC 9197	8/8	99
<i>Penicillium chrysogenum</i>	ATCC 9178	8/8	99



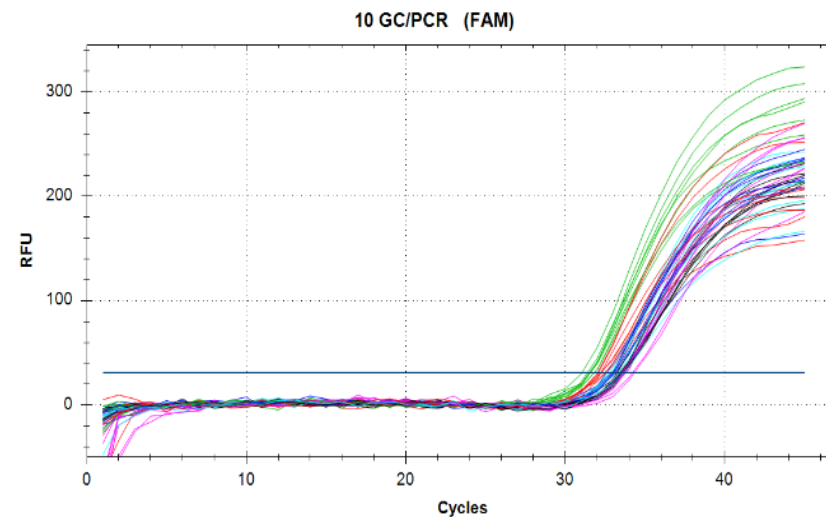
# Limit of detection

Species	Strain No	Acceptance criterion	LOD <sub>95</sub> (CFU/mL)
<i>Bacillus subtilis</i>	ATCC 6633	23/24	25
<i>Clostridium sporogenes</i>	ATCC 19404	23/24	25
<i>Pseudomonas aeruginosa</i>	ATCC 9027	23/24	5
<i>Staphylococcus aureus</i>	ATCC 6538	23/24	10
<i>Streptococcus pyogenes</i>	ATCC 19615	8/8	99
<i>Bacterioides vulgatus</i>	ATCC 8482	8/8	2,5
<i>Escherichia coli</i>	ATCC 8739	8/8	10
<i>Pseudomonas protegens</i>	ATCC 17386	8/8	10
<i>Bacillus cereus</i>	ATCC 10876	8/8	5
<i>Enterococcus faecalis</i>	ATCC 29212	8/8	99
<i>Kocuria rhizophila</i>	ATCC 9341	8/8	10
<i>Staphylococcus epidermidis</i>	ATCC 12228	8/8	99
<i>Serratia marcescens</i>	ATCC 14756	8/8	50
<i>Propionibacterium acnes</i>	ATCC 11827	8/8	25

Species	Strain No	Acceptance criterion	LOD <sub>95</sub> (GC/PCR)
<i>Bacteroides fragilis</i>	ATCC 25285	8/8	10
<i>Enterobacter cloacae</i>	ATCC 13047	8/8	10
<i>Klebsiella pneumoniae</i>	ATCC 13883	8/8	10
<i>Serratia marcescens</i>	ATCC 13880	8/8	10
<i>Clostridium perfringens</i>	ATCC 13124	8/8	10
<i>Yersinia enterocolitica</i>	ATCC 27739	8/8	10

currently tested

recommended extension

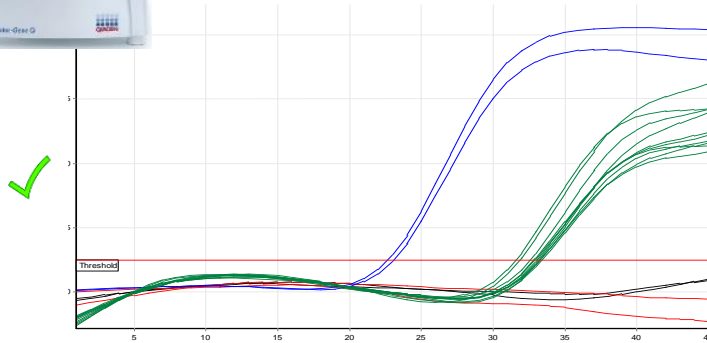




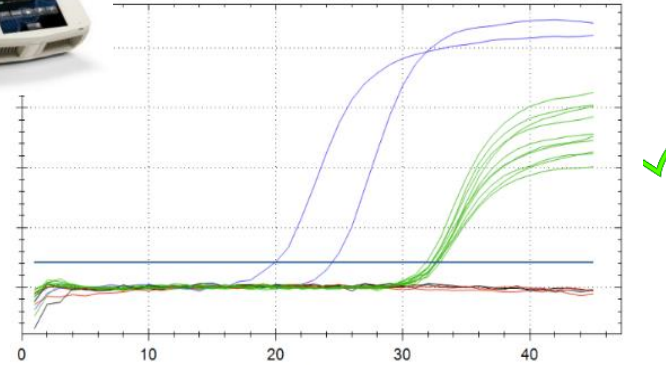
# Device comparability



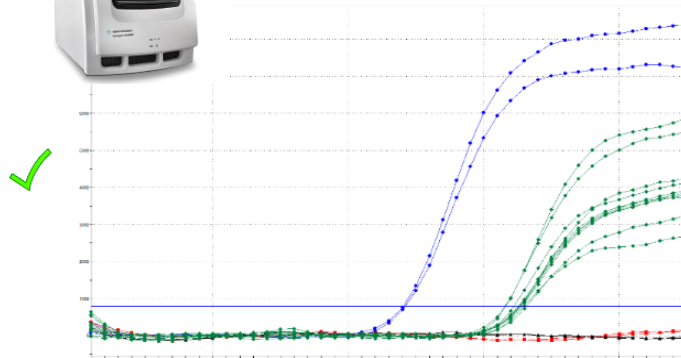
QiaGen Rotor-Gene 6000



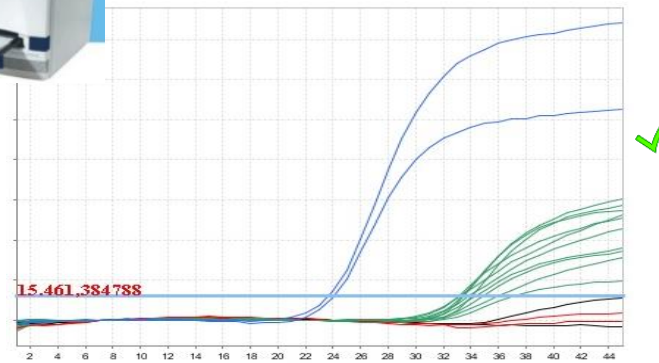
BioRad CFX96 touch



Agilent Mx3005p



Thermo Fischer ABI Prism 7500



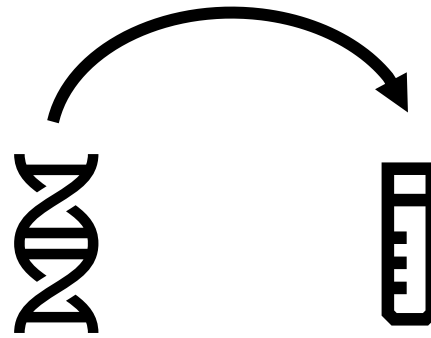
Spiking 99 CFU/ml of the species with the highest LOD<sub>95</sub>  
*Clostridium sporogenes* (LOD<sub>95</sub> = 25)  
*Candida albicans* (LOD<sub>95</sub> = 50)  
Acceptance criterion 8/8 positive

PC  
NTC  
NEC  
Extracts

... by today many more devices are used by our customers.



# False-positives due to free microbial DNA?



*E. coli* DNA spiked into DMEM  
 ↓  
 DNA extraction  
 ↓  
 Real-time PCR

failed		passed							
10 <sup>4</sup> GC/mL		10 <sup>3</sup> GC/mL		100 GC/mL		10 GC/mL		0 GC/mL	
Run 1	Run 2	Run 1	Run 2	Run 1	Run 2	Run 1	Run 2	Run 1	Run 2
39.00	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>
No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>
39.93	39.89	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>
39.38	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>
No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>
No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>
No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>
No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>	No C <sub>q</sub>

There is a low risk of false-positives, because the free DNA is removed efficiently.



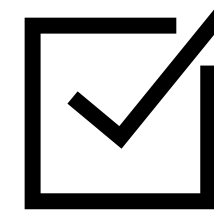
# False-positives due to cell culture medium?

	Bacteria		Fungi	
	Negative Results	% Negative results	Negative Results	% Negative results
DMEM + 5% FCS	8/8	98,95 %	8/8	100 %
	8/8			
	8/8			
	8/8			
	8/8			
	7/8			
	8/8			
	8/8			
	8/8			
	8/8			
	8/8			
	8/8			
	8/8			
DMEM	8/8	100 %	8/8	100 %
	103/104	99 %	104/104	100%

DMEM\* without spike  
acceptance criterion  
> 95 % negative  
\* from Biochrom AG

There is a low risk of false-positives, because cell culture medium is DNA-free.



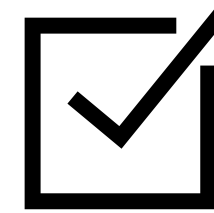


# False-positives due to cell culture medium?

Fungi detection	Negative Results	ROX™	% Negative results
DMEM	8/8	8/8 Correct	100%
Fetal bovine serum (FBS)	8/8	8/8 Correct	100%
DMEM high Glucose, GlutaMAX	8/8	8/8 Correct	100%
DMEM w/o Na-Pyruvat mit stable Glutamin	8/8	8/8 Correct	100%
RPMI 1640	8/8	8/8 Correct	100%
MEM (1)	8/8	8/8 Correct	100%
MEM (2)	8/8	8/8 Correct	100%
KnockOut DMEM	8/8	8/8 Correct	100%
DMEM/F-12 GlutaMAX	8/8	8/8 Correct	100%
Opti-MEM Reduced Serum GlutaMAX	8/8	8/8 Correct	100%
McCoy's 5A Medium	8/8	8/8 Correct	100%
Leibovitz L-15 Medium	8/8	8/8 Correct	100%
Chondrocyte Differentiation Medium	8/8	8/8 Correct	100%
Human Osteoblast Differentiation Medium	8/8	8/8 Correct	100%
ChondroMAX Differentiation Medium	8/8	8/8 Correct	100%
Tahara Lympho One + HS	8/8	8/8 Correct	100%
RPMI, penicillin/Streptomycin, Glutamax, 5% FBS	8/8	8/8 Correct	100%

Cell culture medium  
without spike  
acceptance criterion  
> 95 % negative

There is a low risk of false-positives, because cell culture medium is DNA-free.



## Tolerance of eukaryotic cell background

	<i>Aspergillus brasiliensis</i> 99 CFU (FAM™)			
	$10^6$ cells/ml		$10^5$ cells/ml	
<b>Hela</b>	37.54	31.26	33.72	32.10
	35.24	31.84	33.34	32.07
	<b>4/4</b>		<b>4/4</b>	
<b>Vero</b>	33.30	31.43	32.78	31.39
	34.45	32.76	33.48	31.60
	<b>4/4</b>		<b>4/4</b>	
<b>CHO-K1</b>	33.65	32.40	34.49	33.36
	34.70	32.83	33.89	32.40
	<b>4/4</b>		<b>4/4</b>	
<b>HPBMC</b>	32.21	31.59	33.80	32.12
	32.39	31.25	33.79	31.04
	<b>4/4</b>		<b>4/4</b>	
<b>Jurkat</b>	34.99	34.14	31.77	32.64
	34.80	33.50	34.30	33.81
	<b>4/4</b>		<b>4/4</b>	

All organisms were successfully detected in cell backgrounds of  $10^5$ - $10^6$  cells/ml

# Application Notes

## Rapid, real-time PCR-based detection of microbial contaminations in high cell density Jurkat-, HPBMC-and CHO-cultures using Microsart® ATMP kits

In this study, we assessed the detection capability of Microsart® ATMP Extraction, combined with Microsart® ATMP Bacteria/Fungi/Mycoplasma assays, in high-density cell cultures.



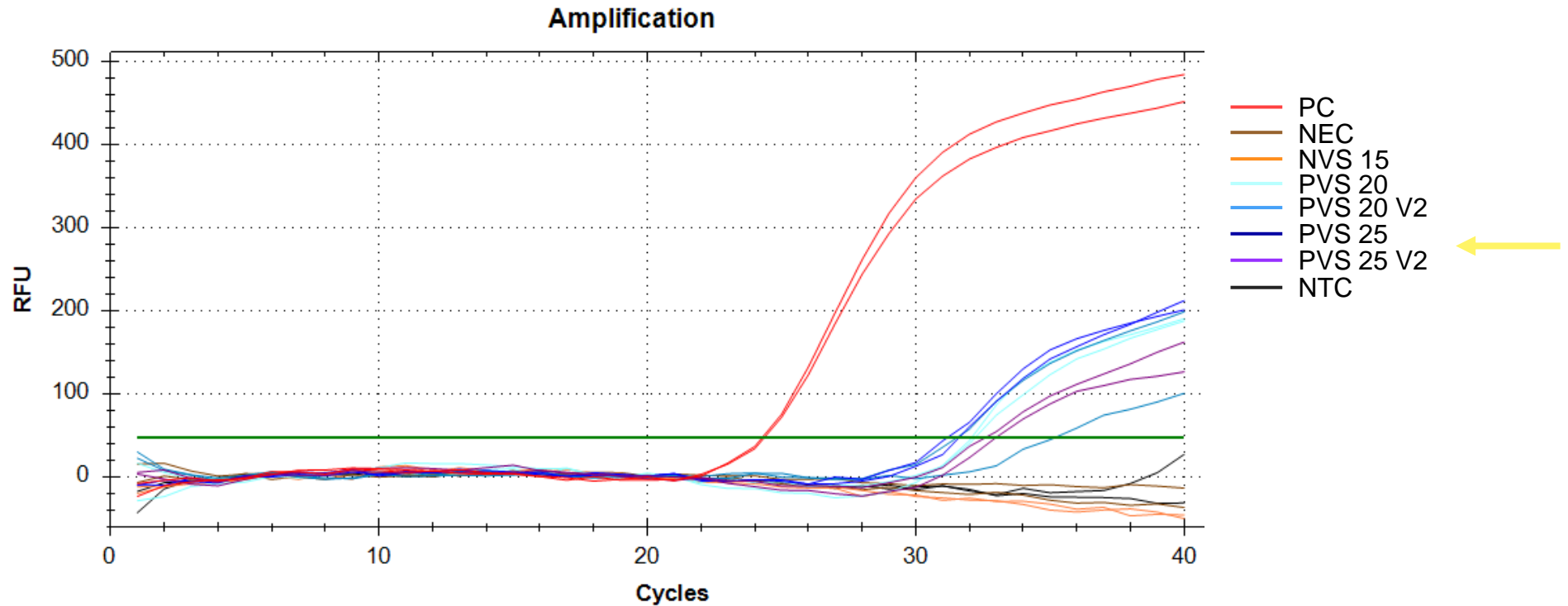
Detection limits in different cell types

Cell Type Background	Microorganism Spike	Background Cells /mL (In 10 <sup>6</sup> )	Detection
CHO	99 CFU <i>B. subtilis</i>	19.0	Successful
CHO	10 CFU <i>M. arginini</i>	15.0 and 15.6 (two individual experiments)	Successful
CHO	10 CFU <i>M. orale</i>	15.5 and 16.3 (two individual experiments)	Successful
Jurkat	99 CFU <i>K. rhizophila</i>	10 to 40	Successful up to 25 x 10 <sup>6</sup> c/mL
Jurkat	50 CFU <i>C. albicans</i>	10 to 40	Successful up to 20 x 10 <sup>6</sup> c/mL
Jurkat	10 CFU <i>M. orale</i>	10 to 40	Not successful: PCR inhibition > 15 x 10 <sup>6</sup> c/mL No detection of Mycoplasma spike
Jurkat	10 CFU <i>M. synoviae</i>	10 to 35	Not successful: Partial PCR inhibition No detection of Mycoplasma spike
HPBMC	99 CFU <i>K. rhizophila</i>	10 to 40	Successful only up to 10 x 10 <sup>6</sup> c/mL
HPBMC	10 CFU <i>M. arginini</i>	15.0	Successful
HPBMC	10 CFU <i>M. orale</i>	19.1	Successful
HPBMC	99 CFU <i>P. aeruginosa</i>	20 and 25	Successful

Table 1: Results of Real-Time PCR detection of respective microbial spikes in varying cell types with respective cell densities.

# Tolerance of eukaryotic cell background

HPBMC cells



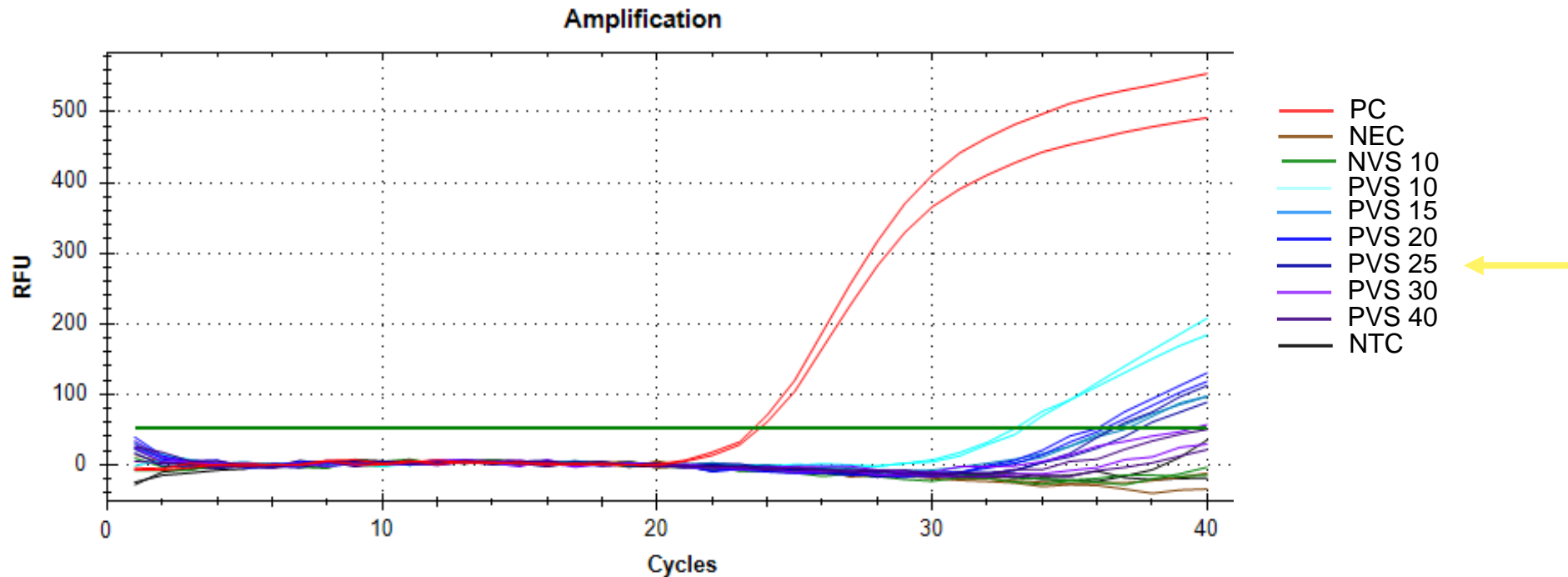
Successful detection of **99 CFU *P. aeruginosa*** in up to **25 million HPBMC** cells.

Medium: Lympho One T-Cell Expansion

HPBMC: human peripheral blood mononuclear cell (e.g. lymphocytes)

# Tolerance of eukaryotic cell background

Jurkat cells



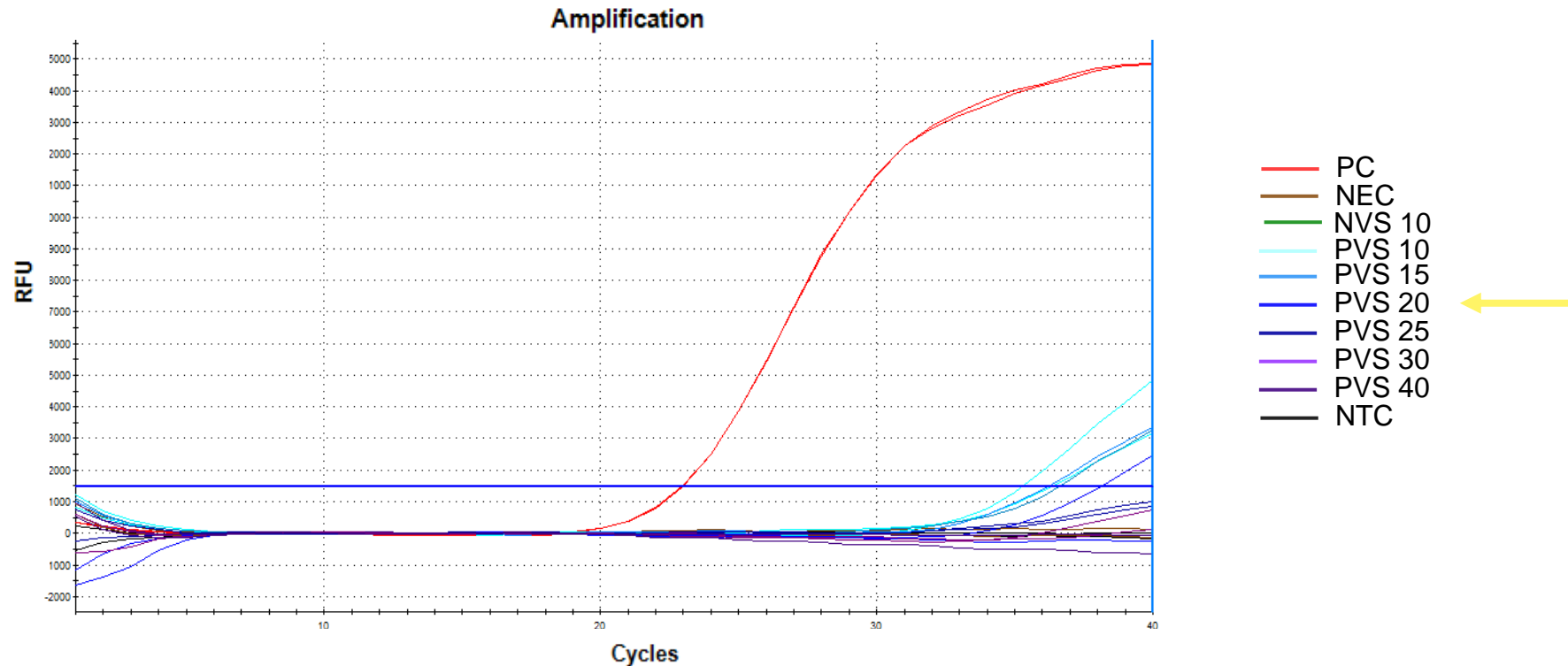
Successful detection of 99 CFU *Kocuria rhizophila* in up to 25 million Jurkat cells.

Medium: RPMI 1640

Jurkat: an immortalized human T lymphocyte cell line

# Tolerance of eukaryotic cell background

Jurkat cells



Successful detection of 50 CFU *C. albicans* in up to 20 million Jurkat cells.

Medium: RPMI 1640

Jurkat: an immortalized human T lymphocyte cell line



# Equivalence with compendial culture method

Sartorius



real-time PCR-based detection

Labor LS

?  
=



classical sterility test according to EP 2.6.1. and USP<71>

2x LOD<sub>95</sub>

### all 6 mandatory species

- *Bacillus subtilis*
- *Clostridium sporogenes*
- *Pseudomonas aeruginosa*
- *Staphylococcus aureus*
- *Candida albicans*
- *Aspergillus brasiliensis*

LOD<sub>95</sub> ½ LOD<sub>95</sub>

### recommended extension

- *Streptococcus pyogenes*
- *Pseudomonas protegens*





# Equivalence with compendial culture method

	Microsart® ATMP Bacteria			Compendial culture method (External)		
	2x LOD <sub>95</sub>	LOD <sub>95</sub>	LOD <sub>95</sub> /2	2x LOD <sub>95</sub>	LOD <sub>95</sub>	LOD <sub>95</sub> /2
<i>Bacillus subtilis</i>	33.16	34.23	35.47	<i>B. subtilis</i>	<i>B. subtilis</i>	<i>B. subtilis</i>
	33.23	34.32	34.38			
<i>Staphylococcus aureus</i>	35.42	35.77	36.56	<i>S. aureus</i>	<i>S. aureus</i>	<i>S. aureus</i>
	34.13	35.67	39.90			
<i>Clostridium sporogenes</i>	34.20	34.87	35.45	<i>C. sporogenes</i>	<i>C. sporogenes</i>	<i>C. sporogenes</i>
	34.10	33.43	35.61			
<i>Pseudomonas aeruginosa</i>	36.40	36.74	37.22	<i>P. aeruginosa</i>	<i>P. aeruginosa</i>	Negative
	36.22	37.96	No Cd			
<i>Streptococcus pyogenes</i>	34.89	35.53	36.55	<i>S. pyogenes</i>	<i>S. pyogenes</i>	<i>S. pyogenes</i>
	35.09	35.93	35.88			
<i>Pseudomonas protegens</i>	34.14	34.38	36.52	Gram - Oxidase +	Gram - Oxidase +	Gram - Oxidase +
	33.28	34.51	35.61			





# Equivalence with compendial culture method

	Microsart® ATMP Fungi			Compendial culture method (External)		
	2x LOD <sub>95</sub>	LOD <sub>95</sub>	LOD <sub>95</sub> /2	2x LOD <sub>95</sub>	LOD <sub>95</sub>	LOD <sub>95</sub> /2
<i>Candida albicans</i>	32.25	32.27	32.96	<i>C. albicans</i>	<i>C. albicans</i>	Negative
	31.94	32.12	33.96			
<i>Aspergillus brasiliensis</i>	34.38	37.06	34.94	<i>A. brasiliensis</i>	<i>A. brasiliensis</i>	<i>A. brasiliensis</i>
	32.40	33.17	34.20			

# Validation overview



- Sensitivity
  - LOD<sub>95</sub> - limit of detection
- Specificity
  - Sequence alignment
  - Sample matrix effects/cross reactivity
  - Specificity of PCR with genomic DNA
  - Comparison with compendial method
- Robustness
  - Spiked cell culture samples
  - Device compatibility
  - Detection of free-DNA
  - False positive rate

Simplifying Progress



Coming  
soon!

Customer validation data  
Service Lab Labor LS

Labor | **LS**

**SARTORIUS**

# RESEARCH

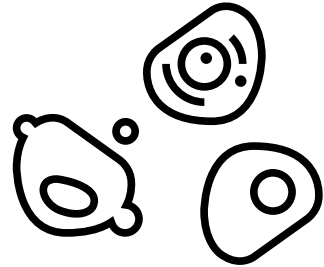


# In-process contamination control of bacteria & fungi

- Key advantages

- Very robust towards inhibitors
- No prior DNA extraction required
- Internal control DNA included in real-time PCR master mix
- One step preparation

→ Quick 'n' Dirty for process monitoring



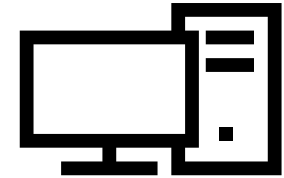
ATMP  
(CHO, HEK,...)



2 µl



23 µl  
RESEARCH  
master mix



Detection in  
real-time PCR cycle:  
Contamination? Yes/No

- Taq-Man® System → reduce false-positive signals
- Duplex assay → reduce false-negative signals
- Universal assay for different real-time PCR cycler → FAM™ and ROX™
- High stability & no freezing → Lyophilized reagents

# In-process contamination control of bacteria & fungi



**Microsart® RESEARCH  
Bacteria**



**Microsart® RESEARCH  
Fungi**

Thank you.

PCR@Sartorius.com



**SARTORIUS**